

## CERTIFICATE OF ANALYSIS

Page 1 of 1

**Product Description:** V-PLEX® and V-PLEX Plus Human IP-10 Kit  
**Kit Lot Number:** K0081773  
**Expiration Date:** 30 Nov 2022  
**Kit Catalog Numbers:** K151NV-Series; K156NV-Series

**Precision and Accuracy Specifications:**

| Spot | Name  | Calibrator |       | Chemokine Panel 1 (human) Control 1 |       | Chemokine Panel 1 (human) Control 2 |       | Chemokine Panel 1 (human) Control 3 |       |
|------|-------|------------|-------|-------------------------------------|-------|-------------------------------------|-------|-------------------------------------|-------|
|      |       | Recovery   | CV    | Recovery                            | CV    | Recovery                            | CV    | Recovery                            | CV    |
| 1    | IP-10 | 80% - 120% | < 20% | 75% - 125%                          | < 20% | 75% - 125%                          | < 20% | 75% - 125%                          | < 20% |

| Spot | Name  | LLOQ  |            |       | ULOQ  |            |       |
|------|-------|-------|------------|-------|-------|------------|-------|
|      |       | Value | Recovery   | CV    | Value | Recovery   | CV    |
| 1    | IP-10 | 1.37  | 80% - 120% | < 20% | 500   | 80% - 120% | < 20% |

**Stock Concentration of Calibrator:**


| Spot | Name  | Conc. | Units |
|------|-------|-------|-------|
| 1    | IP-10 | 2555  | pg/mL |

**Lot-Specific Components:**

| Description                                | Lot Number | Storage Temperature | Expiration Date |
|--|------------|---------------------|-----------------|
| Human IP-10 Plate                          | Z0047346   | 2-8°C               | 30 Nov 2022     |
| Anti-hu IP-10 Antibody                     | D0081401   | 2-8°C               | 30 Apr 2023     |
| Chemokine Panel 1 (human) Calibrator Blend | A0080306   | 2-8°C               | 30 Sep 2026     |
| Diluent 43                                 | M0430045   | ≤ -10°C             | 30 Nov 2023     |
| Diluent 3                                  | M0030109   | ≤ -10°C             | 31 Jan 2024     |

Additional components may be provided to execute the assay. Additional components are not lot-specific.

The above product is intended for Research Use Only. Not for use in Diagnostic Procedures.

|                 | Name          | Function | Signature  | Date        |
|-----------------|---------------|----------|--|-------------|
| Review/Approval | Karen Hamilla | Quality  |  | 23 NOV 2021 |